

Technical Specifications for MicroMounts™

Pins

Length	11, 19, 25 mm and SPINE lengths, marked in 2 mm intervals
Diameter	0.025" (0.64 mm) diameter
Material	Solid 316 non-magnetic stainless steel
Notes	Beveled at 30 degrees at each end.

Crystal holding mount

Material	Polyimide (used in Kapton tape)
Crystal aperture sizes	10, 20, 30, 50, 75, 100, 150, 200, 300, 400, 500 and 600 micrometers
Thickness	7.5, 10 and 12.5 micrometers, depending on size of crystal aperture
Distance from center of alignment cross to center of crystal aperture	870 micrometers
Distance from pin base to center of crystal aperture	12.1 mm (for 11 mm pins) 20.1 mm (for 19 mm pins) 26.1 mm (for 25 mm pins) 18.5 mm (for SPINE length pins)
Width of wicking hole	250 micrometers. Compatible with Size 15 paper wicks.
Length tolerance	0.5 mm for 11, 19 and 25 mm pins, 0.3 mm for SPINE pins.